

Comparison of Whole Blood & Precipitated Blood for the Quantitation of Drugs of Abuse Using PaperSpray

Katherine L. Walker, Yu Zhu, and Neloni Wijeratne

Thermo Fisher Scientific, San Jose, CA 95134

ABSTRACT

Purpose: To compare drug of abuse quantitation in untreated whole blood and whole blood that has been precipitated by zinc sulfate in methanol.

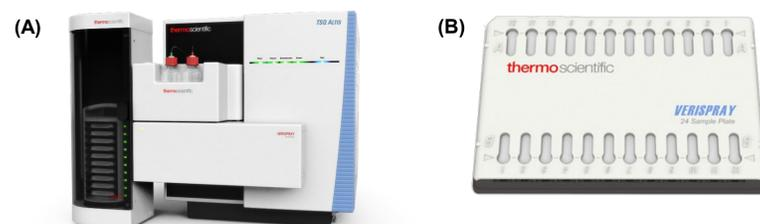
Methods: Twenty-one drugs of abuse and their corresponding internal standards were prepared in whole blood and methanol. Spiked whole blood was mixed in a 1:3 ratio with a precipitation solution (2:1 methanol:0.2 ZnSO₄), centrifuged down, and the supernatant was removed for analysis. The spiked whole blood, precipitated blood, and methanol were deposited onto VeriSpray sample plates, dried, and then analyzed using a Thermo Scientific™ VeriSpray™ PaperSpray ion source coupled to a Thermo Scientific™ TSQ Altis™ MS. The one-minute chromatograms were integrated using TraceFinder 5.1 software and calibration curves were generated.

Results: For drugs in methanol, all had an LLOQ of 5 ng/mL. In whole blood, the signal of calibrators and the background signal dropped due to matrix effects on ionization. For three benzodiazepines, the signal-to-noise and LLOQ was improved by the precipitation method. These compounds either suffer from decomposition or from protein-binding in whole blood. For the remaining 18 drugs of abuse, precipitation does not improve quantitation.

INTRODUCTION

PaperSpray-MS is a technique for rapidly quantifying analytes in dried matrix spots such as whole blood. Little or no sample preparation is required, and sample analysis times are 2 minutes or less. The Thermo Scientific™ VeriSpray™ PaperSpray ion source system utilizes PaperSpray technology to make clinical research workflows faster and more efficient by combining ease-of-use and increased automation with the speed that PaperSpray technology provides. The VeriSpray system consists of the VeriSpray ion source and the Thermo Scientific™ VeriSpray™ plate loader (Figure 1A). Each VeriSpray sample plate contains 24 paper strips (12 on each side, A and B, Figure 1B). The plate loader allows a full 10-plate magazine with a total of 240 samples to be run without user intervention.

Figure 1. (A) VeriSpray ion source and plate loader with loaded magazine mounted to TSQ Altis mass spectrometer (B) VeriSpray sample plate.



Since PaperSpray-MS is a direct analysis technique with no chromatographic separation, the sample matrix can decrease the analyte signal and the LLOQ (lower limit of quantitation). Precipitation of proteins from spiked whole blood with zinc sulfate and extraction with methanol can improve analyte signal, especially for compounds that bind to albumin and other proteins. Furthermore, the matrix effects are lessened from the methanolic solution of precipitated blood compared to whole blood.

In this study we use PaperSpray-MS to analyze 21 drugs of abuse in methanol, whole blood, and precipitated blood to determine the utility of protein precipitation.

MATERIALS AND METHODS

Sample Preparation

- A mix of 21 drugs in the benzodiazepine, opiate, cocaine, stimulant and sedative classes were spiked into human whole blood or methanol (f.c. 5-400 ng/mL) with corresponding internal standards (IS; f.c. 130 ng/mL).
- Samples were put on a blood shaker for 20-30 minutes.
- The precipitated blood sample was prepared from spiked blood: 100 µL of spiked blood was mixed with 300 µL precipitation solution (2:1 methanol: 0.2M ZnSO₄). Vortex and store in fridge for 10 mins. Centrifuge 10 min 12000 rpm. Transfer out 200 µL supernatant.
- For each condition (in methanol, in whole blood, in precipitated blood), 5 replicates of each calibrator level, a matrix blank, and a matrix blank with IS were spotted on VeriSpray sample plates (spotting volume = 8 µL).
- Sample plates were oven-dried at 45 °C for 5 mins and 25 mins for precipitated blood and whole blood, respectively. Samples in methanol dried in 5 mins at r.t.

PaperSpray Conditions

Rewetting (10 µL) and spraying (100 µL) solvents were both 95:5:0.01 methanol: water: acetic acid. The paper tip to MS inlet distance was set to 6.5 mm.

Mass Spectrometer Conditions & Data Analysis

The analysis of drugs of abuse was carried out on a TSQ Altis MS connected with the VeriSpray system. Table 1 and 2 show the MS source parameters and optimized MS transitions, respectively. TraceFinder 5.1 Software was used for processing the 1-minute chromatograms.

Table 1. (A) TSQ Altis MS source conditions and (B) time dependent applied spray voltage setting.

(A) Ion Source Parameter		Value
Spray Voltage	Time Dependent	
Ion Transfer Tube Temp		400 °C
Q1 Resolution		0.7
Q3 Resolution		1.2
CID Gas		1.5 mTorr

(B) Time (min)		Voltage (V)
0		0
0.1		4000
0.95		0

Table 2. Optimized MS transitions for drugs of abuse on TSQ Altis MS with acquisition time of 1.0 min

Compound	Precursor (m/z)	Product (m/z)	Coll. Energy (V)
Norketamine HCl	224.084	124.97	24.55
Norketamine HCl	224.084	179.268	15.49
Ketamine HCl	238.099	124.97	27.87
Ketamine HCl	238.099	207.054	14.35
PCP (Phencyclidine)	244.206	86.125	12.62
PCP (Phencyclidine)	244.206	91.054	31.03
PCP (Phencyclidine)	244.206	159.137	13.97
N-ethylpentylone HCl	250.144	174.125	30.99
N-ethylpentylone HCl	250.144	189.125	23.87
N-ethylpentylone HCl	250.144	202.155	18.61
N-ethylpentylone HCl	250.144	205.125	13.8
Diphenhydramine HCl	256.17	165.125	42.53
Diphenhydramine HCl	256.17	167.125	13.26
Nordiazepam	271.063	140	27.33
Nordiazepam	271.063	208.083	27.33
Diazepam	285.079	153.988	26.52
Diazepam	285.079	193.071	31.03
Diazepam	285.079	222.125	26.15
Diazepam	285.079	257.071	21.76
Hydromorphone	286.144	128.071	55
Hydromorphone	286.144	157.071	38.53
Hydromorphone	286.144	185.071	29.56
Morphine	286.144	152.054	55
Morphine	286.144	165	41.69
Morphine	286.144	181.071	35.79
Morphine	286.144	201.012	24.92
Norhydrocodone HCl	286.144	171.107	36.55
Norhydrocodone HCl	286.144	199.232	26.99
Norhydrocodone HCl	286.144	241.107	23.7
Benzoylcegonine	290.139	105.125	29.14
Benzoylcegonine	290.139	168.196	18.73
Hydrocodone	300.159	128.143	55
Hydrocodone	300.159	171.125	40.17
Hydrocodone	300.159	199.125	29.18
Hydrocodone	300.159	243.083	22.9
Codeine	300.159	165.071	39.54
Codeine	300.159	215.012	25.85
Codeine	300.159	225.083	27.41
Temazepam	301.074	176.97	37.98
Temazepam	301.074	193.054	33.85
Temazepam	301.074	239.042	46.95
Temazepam	301.074	255.054	21.64
Oxymorphone	302.139	198.125	41.77
Oxymorphone	302.139	227.083	27.75
Cocaine	304.154	82.125	29.47
Cocaine	304.154	182.179	18.94
Zolpidem	308.176	219.095	53.82
Zolpidem	308.176	221.155	38.53
Zolpidem	308.176	235.137	34.15
Alprazolam	309.09	205	40.59
Alprazolam	309.09	274.226	24.33
Alprazolam	309.09	281.071	25.35
Clonazepam	316.048	190.202	42.74
Clonazepam	316.048	214.071	36.09
Clonazepam	316.048	269.982	24.42
Midazolam	326.085	222.083	46.91
Midazolam	326.085	249	36.97
Midazolam	326.085	291.083	26.19
Zolpidem Phenyl-4-carboxylic acid	338.15	219.137	54.24
Zolpidem Phenyl-4-carboxylic acid	338.15	265.083	35.71
Zolpidem Phenyl-4-carboxylic acid	338.15	293	26.19

RESULTS

Calibration curves for each drug of abuse were generated for analyte in methanol, blood, and precipitated blood. The lower limit of quantification (LLOQ) was set to the lowest calibration standard analyzed that yielded < 20% accuracy and < 15% CV for 5 replicate samples. Example calibration curves of a representative drug in the benzodiazepine, opiate, cocaine, and sedative classes in whole and precipitated blood are shown in Figure 2.

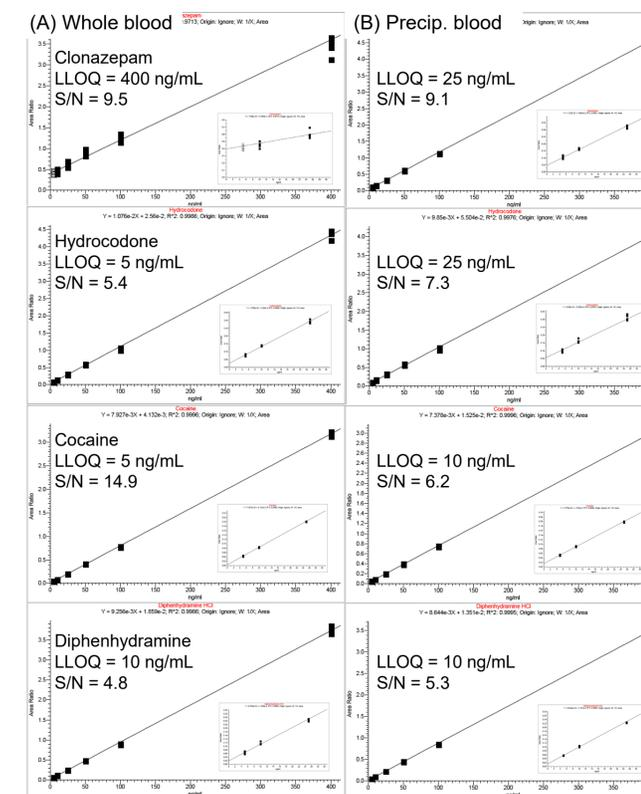
Table 3. LLOQs for drugs of abuse in methanol (neat), in blood, and in precipitated blood.

Compound	LLOQ Neat (ng/mL)	LLOQ in Whole Blood (ng/mL)	LLOQ in Precip Blood (ng/mL)
Alprazolam	5	10	5
Benzoylcegonine	5	10	25
Clonazepam	5	400	25
Cocaine	5	5	10
Codeine	5	25	100
Diazepam	5	25	25
Diphenhydramine	5	10	10
Hydrocodone	5	5	25
Hydromorphone	5	50	50
Ketamine	5	5	25
Midazolam	5	5	5
Morphine	5	25	50
N-ethylpentylone	5	25	25
Nordiazepam	5	5	5
Norhydrocodone	5	10	50
Norketamine	5	10	25
Oxymorphone	5	25	50
PCP	5	5	5
Temazepam	5	400	25
Zolpidem Phenyl-4-carboxylic acid	5	5	5
Zolpidem	5	5	5

In methanol, the LLOQ was 5 ng/mL, the lowest level calibration standard. In whole blood, the signal of calibrators and the signal of the background decreased compared to results in methanol due to matrix effects. The signal-to-noise was reduced for analytes in blood. In precipitated blood, the signal of the background was between the values for those in whole blood and methanol. Depending on the analyte, the signal of calibrators either increased or remained the same when compared to the signal in whole blood. For three benzodiazepines—alprazolam, clonazepam, and temazepam—the LLOQ in precipitated blood decreased and the linearity of the calibration curve improved when compared to whole blood (Table 3). For nine analytes—mostly benzodiazepines, stimulants and sedatives—the LLOQ remained the same for whole blood and precipitated blood. For the remaining nine analytes—mostly opiates and cocaine—the LLOQ decreased up to 4-times due to the 4-fold dilution during the preparation of analyte in precipitated blood.

RESULTS CONT.

Figure 2. Calibration curves of clonazepam, hydrocodone, cocaine, and diphenhydramine in (A) whole blood and (B) precipitated blood. Inset: zoom from 1-30 ng/mL.



CONCLUSIONS

- PaperSpray is a sensitive, direct technique for the analysis of drugs of abuse in whole blood. Sample preparation is not necessary for a majority of analytes.
- For benzodiazepines, precipitation with zinc sulfate and methanol is a simple, rapid clean-up method to improve the LLOQ. In particular, clonazepam and temazepam improved significantly in precipitated blood because they bind strongly to albumin.
- For difficult compounds, especially known protein-binders, precipitation with zinc sulfate is a simple, rapid clean-up method that should be considered when optimizing sample preparation.

REFERENCES

- Espy, R.D. et al. Paper spray and extraction spray mass spectrometry for the direct and simultaneous quantification of eight drugs of abuse in whole blood. *Anal Chem.* **2014**, *86*, 15, 7712-7718.
- Chin, P.K.L. et al. Adult age and *ex vivo* protein binding of lorazepam, oxazepam and temazepam in healthy subjects. *Br J Clin Pharmacol.* **2011**, *72*, 6, 985-989.
- Pacifici, G.M. et al. Plasma protein binding of clonazepam in hepatic and renal insufficiency and after hemodialysis. *Ther Drug Monit.* **1987**, *9*, 4, 369-373.

TRADEMARKS/LICENSING

© 2020 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others.

PO65747-EN 0422S

ThermoFisher
SCIENTIFIC